



# SPECIFICATIONS

## CREOPTIX WAVE

### GENERAL

Noise (RMS)	<0.01 pg/mm <sup>2</sup> @ 1 Hz
Drift	<0.3 pg/mm <sup>2</sup> /min
Readout Frequency	1 Hz, 10 Hz or 40 Hz
Association Const. Range	$k_a = 10^3 - 5 \times 10^7 \text{ M}^{-1} \text{ s}^{-1}$ (small molecules) $k_a = 10^3 - 3 \times 10^9 \text{ M}^{-1} \text{ s}^{-1}$ (large molecules)
Dissociation Const. Range	$k_d = 10^{-5} - 10 \text{ s}^{-1}$
Analysis Temperature Range	15°C - 40°C
Molecular Weight Limit	No lower limit

waveRAPID® Functionality No

### FLUIDICS

Flow Channels / Path	2, parallel
Channel Referencing	2-1 and 1-2
Flow Cells	Sealed, disposable, integrated into disposable WAVEchip
Flow Rate	1 - 400 µl/min
Crude Sample Robustness	Yes

### SAMPLE HANDLING

Sample Capacity	2x microtiter plates (96 or 384 well, standard or deep well) or vial racks (48 positions of 1.5ml)
Buffer	1 buffer
Degasser	Built-in
Injection Volume	< 450 µl, 100 µl typical
Sample Volume Required	Injection volume plus 15-50 µl (application dependent)
Sample Storage Temperature	Ambient or 4°C - 20°C regulated
Sample Recovery	Yes
Automation	120h of unattended operation

### DATA TREATMENT

Information Provided	Kinetic and affinity data ( $k_a$ , $k_d$ , $K_D$ )
Graphs	Real-time curves, multiple curve overlays, fit, report point plots
Data Extraction	Curves, $k_a$ , $k_d$ , $K_D$ tables, graphs, reports
Data Analysis	Fully automated data evaluation
Kinetic Models	Predefined models including 1:1 interaction, mass transport, heterogenous ligand, conformational change and bivalent
Direct Kinetics	Yes

